

Histone H3T32ph antibody (mAb)

Catalog Nos: 61777, 61778

RRID: AB_2793762

Clone: 6C7G12

Isotype: IgG2a

Application(s): ICC, IF, WB

Reactivity: Human, Wide Range Predicted

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 17 kDa

Background: Histone H3 is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 147 base pairs of DNA wrapped around an octamer of core histone proteins (two each of Histone H2A, Histone H2B, Histone H3 and Histone H4). Chromatin is subject to a variety of chemical modifications, including post-translational modifications of the histone proteins and the methylation of cytosine residues in the DNA. Reported histone modifications include acetylation, methylation, phosphorylation, ubiquitylation, glycosylation, ADP-ribosylation, carbonylation and SUMOylation; these modifications play a major role in regulating gene expression.

Phosphorylation of several residues of histone H3, such as Ser10 (Histone H3 phospho Ser10), Ser28 (Histone H3 phospho Ser28) and Thr11, is tightly correlated with chromosome condensation during both mitosis and meiosis. Phosphorylation of Thr32 has been reported in both plants and animals and is enhanced in cells undergoing mitosis. However, its distribution on the chromosome appears different than these other H3 phosphorylation events suggesting phosphorylation of Thr32 may have a distinct role.

Immunogen: This antibody was raised against a peptide containing phosphothreonine 32 in human Histone H3.1.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

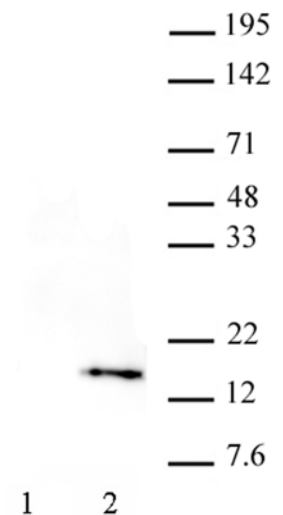
Application Notes:

Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Histone H3T32ph antibody (mAb) tested by Western blot.

HeLa acid extract (10 µg per lane) was probed with Histone H3T32ph antibody at a dilution of 1 µg/ml.

Lane 1: No treatment.

Lane 2: Cells treated with colcemid.